

Certificate of Analysis

Kit Name:	Simoa™ Aβ42 2.0 Kit
Product Number:	101664
Lot Number:	500823
Expiration:	19-Jan-2018

113 Hartwell Ave.
Lexington MA
02421

www.quoterix.com

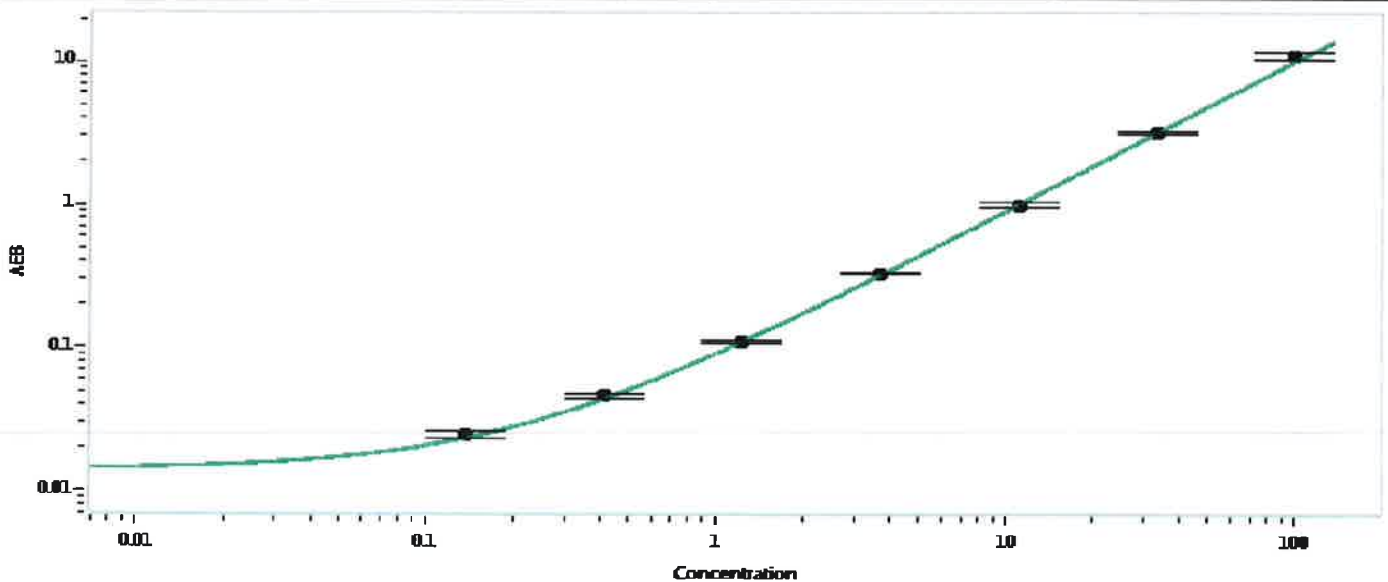
Component	Part Number	Lot Number
Aβ 42 2.0 Bead Reagent	101666	71243
Aβ 42 2.0 Detector Reagent	101665	71114
Aβ 42 2.0 SBG Reagent	101667	71113
Aβ 42 2.0 Sample Diluent	101668	71092
Aβ 42 Calibrator Concentrate	100717	71660
Aβ 42 2.0 Calibrator Diluent	101669	71093
RGP Reagent	100030	N/A ¹

¹ RGP is not kit lot specific

Calibrator Concentrate Value Assignment (pg/mL)	
Assigned Value	1991

Controls	Lot Number	Range (pg/mL)	Value at Release (pg/mL)
Control 1	52520	0.681-1.39	1.38
Control 2	52520	27.1-40.6	34.3

Patient Panels	Lot Number	Range (pg/mL)	Value at Release (pg/mL)
Panel 1	MF-082015-01	0.942-1.53	1.07
Panel 2	MF-081415-02	7.46-11.2	8.42
Panel 3	MF-081415-03	9.20-15.0	10.6
Panel 4	MF-090315-09	27.6-66.8	45.8



Review/Approval

David Rice

QC Manager

DDR 26 Jun 2017

Name

Title

Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications. It is intended for Research Use Only. Not for use in Diagnostic Procedures.